

---

# PAK4

Product code: 07-126

Full-length human PAK4 [1-591(end) amino acids of accession number NP\_005875.1] was expressed as N-terminal GST-fusion protein (91 kDa) using baculovirus expression system.

Assay platform : MSA  
Substrate : SGKtide  
  
ATP ( $\mu\text{M}$ )  $K_m$  app /  $B_{in}$  : 2.5 / 5  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP  $B_{in}$  (nM) : 12  
IC50 at 1 mM ATP (nM) : n.a.

---